

Instruction Manual

Catalog Number	PK-CA577-1042-1G
Description	An anti-inflammatory glucocorticoid. Inhibits the expression of the inducible but not the constitutive nitric oxide synthase in vascular endothelial cells ($IC_{50} = 5 \text{ nM}$). Enhances active cation transport in aortic smooth muscle cells by stimulating the Na^+K^+ pump. Has antiinflammatory and anti-rheumatic properties. Induces apoptosis in human thymocytes. In general, 500-1000 nM of dexamethasone is sufficient to induce apoptosis following a 6-hour incubation at 37°C.
Quantity	1 g
Sequence / Molecular Weight / Molecular Formula	392.5 Da; $C_{22}H_{29}FO_5$
Purity	>99% by TLC
Appearance / Formulation / Solubility	White solid (m.p. 262-264°C). Soluble in ethanol (25 mg/ml).
Storage & Stability	Store powder at desiccated at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Applications	see Description
References	1. Fearnhead, H.O., et al. (1994) <i>Biochem. Pharmacol.</i> 48:1073-1079. 2. Moncada, S., et al. (1991) <i>Trends Pharmacol. Sci.</i> 12:130-143.
Caution	NA

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.